

## Nicotinic ACh7R activation and pharmacology recorded on Nanion's Patchliner

The electrophysiology team at Nanion Technologies GmbH, Munich. Cells kindly provided by Galantos Pharma, GmbH, Germany.



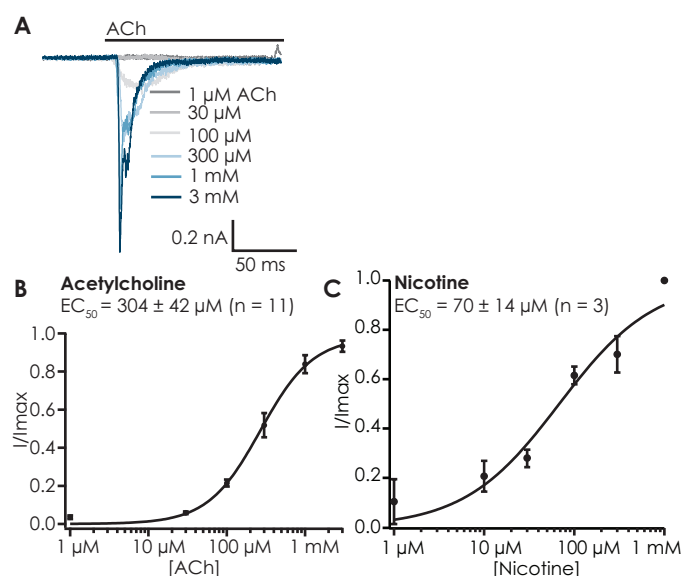
### Summary

The nicotinic acetylcholine receptor (nAChR) is a member of the ligand-gated ion channel superfamily which includes GABA<sub>A</sub>, 5HT<sub>3</sub>, NMDA and glycine receptors. It is a cation-permeable ion channel activated by the neurotransmitter acetylcholine and the natural alkaloid, nicotine. Neuronal nAChR are pentameric and functional channels are formed from a repertoire of nine  $\alpha$  ( $\alpha 2$  to  $\alpha 10$ ) and three  $\beta$  subunits ( $\beta 2$  to  $\beta 4$ ). Most nAChR exist as heteromers with the stoichiometry 2 $\alpha$  to 3 $\beta$ , however some  $\alpha$  subunits function as homomers, these being  $\alpha 7$  or  $\alpha 9$  (for reviews see Refs. 1 & 2). nAChR have been proposed to play a role in many neurological disorders such as Alzheimer's Disease, Parkinson's, schizophrenia and depression<sup>1-4</sup>. nACh7R are widely distributed in the mammalian brain including in the cerebral cortex, hippocampus, basal ganglia and cerebellum<sup>4</sup>. There is evidence that nACh7R play a role in cognition<sup>5,6</sup> and could be a potential therapeutic target in cognitive disorders such as Alzheimer's Disease or schizophrenia<sup>3,5,6</sup>.

Here we present data collected on a Patchliner showing the potential use of the Patchliner to record fast desensitizing nACh7R currents activated by acetylcholine (ACh) or nicotine. ACh and nicotine activated nACh7R in a concentration-dependent manner with an EC<sub>50</sub> similar to those reported in the literature<sup>7-10</sup>. nACh7R could be repetitively activated by nicotine and were enhanced by the Type I positive allosteric modulator (PAM), NS1738, with an EC<sub>50</sub> in good agreement with the literature<sup>11</sup>.

### Results

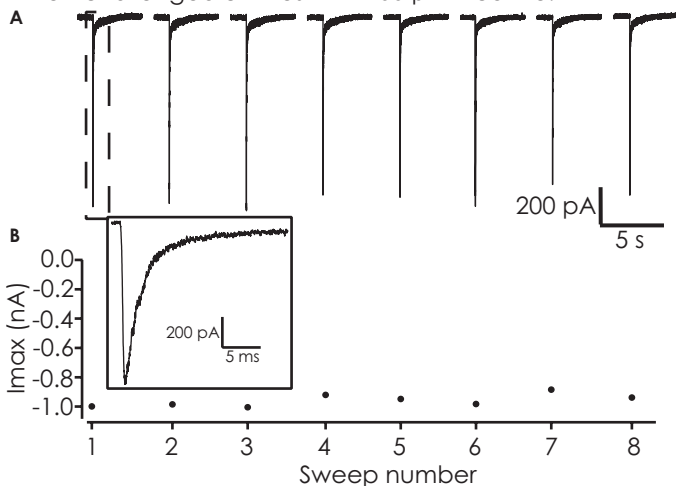
Nicotinic ACh7R were activated by increasing concentrations of ACh or nicotine (Figure 1). A concentration response curve for ACh (Fig. 1B) revealed an EC<sub>50</sub> of 304 ± 42  $\mu$ M (n = 11) in good agreement with the range found in the literature<sup>8-10</sup>. Raw traces from an example cell are shown in Fig. 1A. The concentration response curve for nicotine is shown in Fig. 1C which reveals an EC<sub>50</sub> = 70 ± 14  $\mu$ M (n = 3), in good agreement with the literature<sup>7-9</sup>.



**Figure 1:** Activation of nACh7R by increasing concentrations of ACh or nicotine. **A** Raw traces from an example cell showing activation of nACh7R by increasing concentrations of ACh. **B** Concentration response curve for ACh activation for an average of 11 cells, EC<sub>50</sub> = 304 ± 42  $\mu$ M (n = 11). **C** Average concentration response curve for nicotine activation, EC<sub>50</sub> = 70 ± 14  $\mu$ M (n = 3).

# Application Note

Figure 2 shows the repetitive activation of nAChA7R. Currents were activated with a similar peak amplitude when challenged 8 times with 100  $\mu\text{M}$  nicotine.



**Figure 2:** **A** Nicotinic AChA7R could be repetitively activated by 100  $\mu\text{M}$  nicotine. Inset shows nAChA7R activation of the first application expanded **B** Timecourse of the experiment. Nicotinic AChA7R currents of approx. 1 nA were activated 8 times reproducibly in the same cell.

NS1738 is a Type I PAM of nAChA7R which increases amplitude without any profound effects on desensitization<sup>11</sup>. It has been shown to produce cognitive enhancement in rats. We show that NS1738 enhanced nAChA7R-mediated responses when pre-

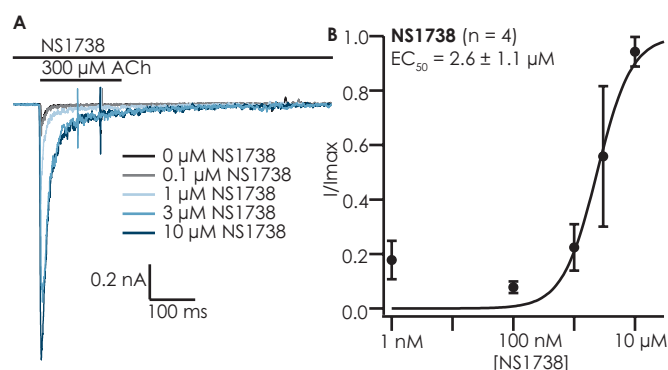
## References

1. Colquhoun *et al.*, in Burger's Medicinal Chemistry, Vol II Drug Discovery and Drug Development, Chapter 11, pp 357-405. Ed. Abraham, D., J. Wiley, New York.
2. Albuquerque *et al.*, 2009. *Physiol Rev.* 89: 73–120
3. Araki *et al.*, 2002. *Jpn. J. Pharmacol.* 88: 133 – 138
4. Hurst *et al.*, 2013. *Pharmacol. & Therapeut.* 137: 22–54
5. Leiser *et al.*, 2009. *Pharmacol. & Therapeut.* 122: 302–311
6. Pohanka, M. 2012. *Int. J. Mol. Sci.* 13: 2219-2238
7. Fu *et al.*, 2008. *Am. J. Resp. Cell & Mol. Biol.* 41: 94-99
8. Corradi & Bouzat. 2016. *Mol. Pharmacol.* 90:288–299
9. Virginio *et al.*, 2002. *Eur. J. Pharmacol.* 445: 153–161
10. Andersen *et al.*, 2013. *PNAS.* 110 (51): 20819–20824
11. Timmermann *et al.*, 2007. *JPET* 323: 294–307

## Methods

### Cells

An inducible HEK293 cells stably expressing human nAChA7R were used.



**Figure 3:** Enhancement of nAChA7R by NS1738. NS1738 at increasing concentrations (0.1  $\mu\text{M}$  - 10  $\mu\text{M}$ ) was pre-incubated and then co-applied with 300  $\mu\text{M}$  ACh. NS1738 enhanced the peak of the ACh response at a concentration of 1  $\mu\text{M}$  or above. **B** Concentration response curve for NS1738 enhancement,  $EC_{50} = 2.6 \pm 1.1 \mu\text{M}$  ( $n = 4$ ).

incubated and then co-applied with acetylcholine (Fig. 3). A full concentration response curve to NS1738 was performed revealing an  $EC_{50}$  of  $2.6 \pm 1.1 \mu\text{M}$  ( $n = 4$ ), in excellent agreement with the literature<sup>11</sup>.

In summary, nAChA7R stably expressed in HEK293 cells can be reliably activated by nicotine or ACh and enhanced by the Type I PAM, NS1738, with  $EC_{50}$  values in good agreement with published literature<sup>7-11</sup>. Therefore, the Patchliner provides a viable, higher throughput alternative to conventional patch clamp for the discovery of agonists of nAChA7R as potential therapeutics to enhance cognitive function.

### Cell culture

Cells were cultured and harvested according to Nanion's standard cell culture protocol.

### Electrophysiology

Whole cell patch clamp recordings were conducted according to Nanion's standard procedure for the Patchliner. Cells were held at a holding potential of -80 mV. To achieve short exposure times, solutions were stacked in the robotic pipettor. First, wash solution (280  $\mu\text{l}$ ) was aspirated followed by aspiration of the agonist-containing solution (30  $\mu\text{l}$ ) and then rapidly applied to the cell at a speed of 171  $\mu\text{l/s}$ . Wash solutions contained the PAM in the NS1738 experiments. The cells were pre-incubated in NS1738 before co-application with 300  $\mu\text{M}$  ACh.